

Datasheet: LNK262B

Description:	LYNX RAPID PLUS BIOTIN (TYPE 1) ANTIBODY CONJUGATION KIT
Name:	BIOTIN (TYPE 1) CONJUGATION KIT
Format:	Kit
Product Type:	Conjugation Kit
Quantity:	3 CONJUGATIONS for 200µg antibody

Product Details

Applications

This product has been reported to work in the following applications. This information is derived from testing within our laboratories, peer-reviewed publications or personal communications from the originators. Please refer to references indicated for further information. For general protocol recommendations, please visit www.bio-rad-antibodies.com/protocols.

	Yes	No	Not Determined	Suggested Dilution
Conjugation	▪			

We recommend that for each conjugation the user determines the best antibody:conjugate ratio.

Product Information

LYNX Rapid Plus Biotin (Type 1) Antibody Conjugation Kit® enables the rapid conjugation of a pre-prepared lyophilized mixture containing Biotin label to an antibody or protein in minutes. Activation of proprietary reagents within the antibody-label solution results in the coupling of the antibody to Biotin.

The LYNX Rapid Plus Conjugation kit® can be used to label small quantities of antibody/protein at near neutral pH, allowing a high conjugation efficiency with 100% antibody recovery and no requirement for desalting or dialysis.

This kit has been optimized for the additional use of a streptavidin labelled reagent. For customers intending to capture their conjugate on Streptavidin coated plates, we recommend our [LYNX Rapid Plus Biotin \(Type 2\) Antibody Conjugation Kit](#) instead.

Reagents In The Kit

3 Vials of 100ug LYNX Rapid Plus lyophilized Biotin mix
1 Vial LYNX Modifier reagent
1 Vial LYNX Quencher reagent

Preparing The Antibody

The following buffer solutions are recommended:for preparing the antibody:

10-50mM amine-free buffer (e.g HEPES, MES, MOPS and phosphate) pH range 6.5-8.5, although moderate concentrations of Tris buffer (<20mM) may be tolerated.

Do not use buffers containing nucleophilic components (e.g primary amines), thiols (e.g. Thiomersal/Thimerosal), Merthiolate, Glycine or Proclin. These substances may react with LYNX chemicals. Azide (0.02-0.1%), EDTA and common non-buffering salts and sugars have little or no effect on conjugation efficiency.

It is recommended that up to 100-200 µg antibody be used in each labeling reaction. For optimal results the antibody volume should be 40-100 µl, at a concentration range between 1mg/ml and 2.5

mg/ml.

-
- Instructions For Use**
1. Before you add antibody to the Lynx Rapid Plus mix, add 1 µl of Rapid Modifier reagent for each 10 µl of antibody to be labelled. Mix gently.
 2. Remove the screw cap from the vial of Lynx Rapid Plus mix and pipette the antibody sample (with the added Rapid Modifier) directly onto the lyophilised material. Resuspend gently by withdrawing and re-dispensing the liquid once or twice using a pipette.
 3. Replace cap on the vial and leave standing for 15 minutes at room temperature (20-25 °C). Longer incubation times could have negative effects on the conjugate.
 4. After 15 minutes incubation, add 1 µl of Lynx Rapid Plus Quencher reagent for every 10 µl of antibody used. The conjugate can be used after 4 minutes. The conjugates do not require purification.

Storage

Store kit at -20°C only.
Newly-conjugated antibody can be stored at 4°C. For long term storage however, the addition of a preservative is recommended.
Storage in frost-free freezers is not recommended.
This product should be stored undiluted.
Avoid repeated freezing and thawing.

Shelf Life 12 months after despatch.

Health And Safety Information

Material Safety Datasheet documentation #10533 #10567 #10568 available at:
Lyophilized Biotin Mix (Type 1) (10533): <https://www.bio-rad-antibodies.com/uploads/MSDS/10533.pdf>
Rapid Plus Quencher Reagent (10567): <https://www.bio-rad-antibodies.com/uploads/MSDS/10567.pdf>
Rapid Plus Modifier Reagent (10568): <https://www.bio-rad-antibodies.com/uploads/MSDS/10568.pdf>

Licensed Use

Bio-Rad LYNX Rapid Plus Conjugation kits® are offered for research purposes alone, and are not intended for human, therapeutic or diagnostic use. The purchase of this conjugation kit conveys to the buyer (whether the buyer is a not-for-profit, academic or for-profit entity) the non-transferable right to use the amount of product purchased and the components of the product for in-house research. The buyer shall not sell or otherwise transfer this product, its components, or materials prepared therefrom to any third party. The buyer shall not use this product or its components for commercial purposes. For the avoidance of doubt, 'commercial purposes' means any activity by a party for consideration and includes, without limitation, use of the product or its components (i) in the manufacturing of conjugated materials (e.g. labeled antibodies), (ii) to provide a service, information or data, (iii) for therapeutic, diagnostic or prophylactic purposes, or (iv) for repackaging/resale, whether or not such product or its components are resold for use in research. The use of this product by the buyer constitutes agreement with the terms of this limited use label license for LYNX products.

Regulatory For research purposes only

North & South America Tel: +1 800 265 7376

Fax: +1 919 878 3751

Email: antibody_sales_us@bio-rad.com

Worldwide

Tel: +44 (0)1865 852 700

Fax: +44 (0)1865 852 739

Email: antibody_sales_uk@bio-rad.com

Europe

Tel: +49 (0) 89 8090 95 21

Fax: +49 (0) 89 8090 95 50

Email: antibody_sales_de@bio-rad.com

'M317653:180705'

